



sartorius

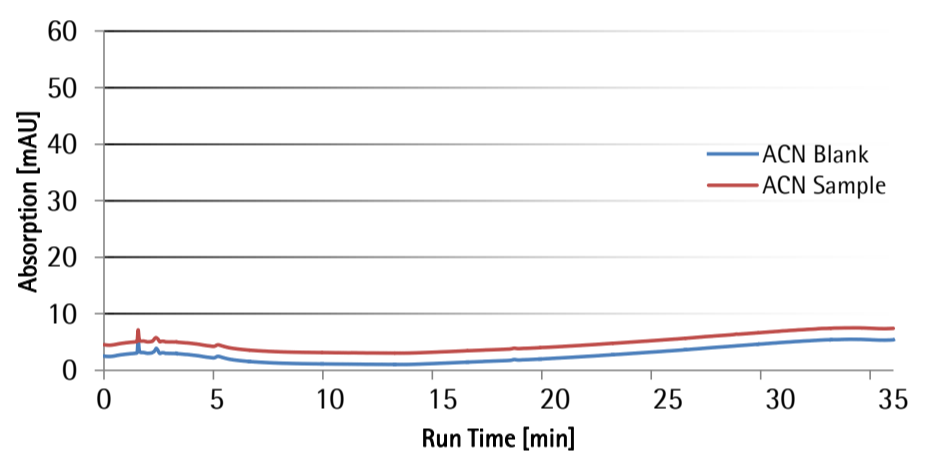
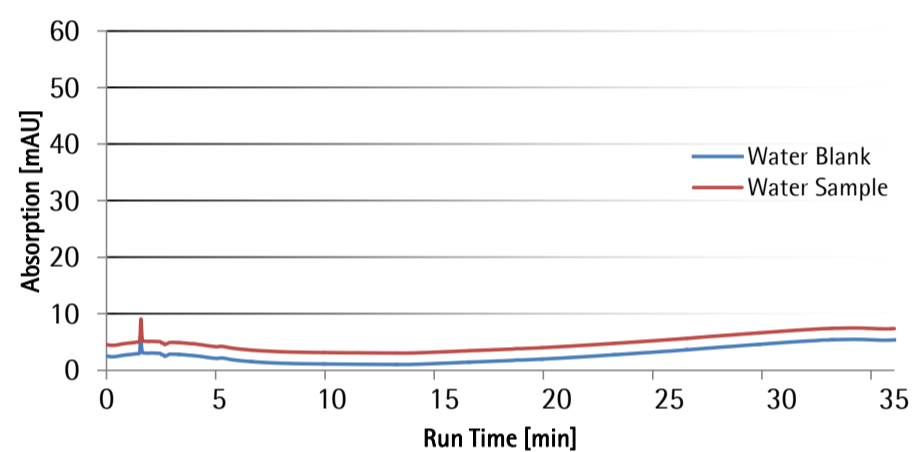
HPLC Purity Certificate for Minisart[®] RC

Order Number: 17765 ----- Q

Lot Number: 00683103

This is to certify that the product described above has been analyzed for extractables by testing the incorporated membrane by HPLC-UV/Vis.

Chromatograms of the HPLC-UV/Vis analyses of the extracts



HPLC Method

Column: C18, 250 x 4.6 mm, Column Temperature: 40°C, Injection Volume: 20 µL, Flow Rate: 1 mL/min,
Detection Wavelength: 220 nm, Analysis Time: 35 min

Mobile Phases: A) Acetonitrile, B) Water

Gradient: 60% A for 10 min, 60% to 100% A in 20 min, 100% A for 5 min

Filter device eluted with 0.5 mL extraction solvent per cm² membrane area (Surface-Area to Volume Ratio: 2 cm²/mL)

For regenerated cellulose (RC), the extraction strength of the solvent acetonitrile has been tested to be comparable to methanol.

2020-06-22



Dr. Anna Vreemann
Site Quality Manager

Manufactured by Sartorius Stedim Biotech GmbH
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289